

Europe Office and Corporate Headquarters  
Cisbio Bioassays  
Phone: +33 (0)4 66 79 67 05  
Fax: +33 (0)4 66 79 19 20  
E-mail: bioassays@cisbio.com

Japan Office  
Sceti Medical Labo K.K.  
Phone: +81 (0)3 5510 2932  
Fax: +81 (0)3 5510 0130  
E-mail: reagent@scetimedilabo.co.jp

USA Office  
Cisbio US, Inc.  
Phone : 888 963 4567  
Fax : 781 687 1500  
E-mail : htrfinfo@cisbio.us



## Tag-lite GIP receptor red ago. - 5,000 tests

**Product number:** L0018RED  
**Batch number :** 1  
**Storage temperature:** -20°C

[www.htrf.com](http://www.htrf.com)

## Product Information

For in vitro research use only

Version Mar. 11 - D

<b>Formulation</b>	100 µL, 10 µM solution, H2O
<b>Size</b>	5,000 tests* - GIPR assay
<b>MW</b>	5851.4
<b>Measure</b>	665 nm

\* 5,000 tests in a competition assay at the recommended concentration in a 384-well format (20 µL low volume)

## Description

This agonist is a GIP derivative labeled with a Red Emitting HTRF fluorescent probe.

## Specificity

This ligand has been successfully used in a GIP receptor Tag-lite binding assay.

## Characterization

Receptor	Assay format	Kd	Working concentration*
GIPR	HEK293T / 384SV suspension	14 +/- 5 nM	10 nM

\* Working concentration advised for ligand competition assays.

## Storage and stability

Upon receiving the ligand, store at -20°C or colder and away from bright light.  
Avoid repeated freeze-thaw cycles, as this can degrade the ligand. Individual aliquots should be prepared.

## Protocole

Information available on: [www.htrf.com](http://www.htrf.com)

## MSDS

Information available on: [www.htrf.com](http://www.htrf.com)

Copyright © 2010 Cisbio Bioassays, France

Lumi4 is a registered trademark of Lumiphore, Inc. Lumi4-based products may be covered under one or more of U.S. patents: 6,515,113, 6,864,103, 7,442,558, 6,406,297, 7,018,850, and 7,404,912 or international counterparts.

HTRF®, Tag-lite®, and the HTRF® logo are trademarks belonging to Cisbio Bioassays.

HTRF® products are manufactured under one or more of the following patents and foreign equivalent: EP 0 180 492 / US 4,927,923 / US 5,220,012 / US 5,432,101 - EP 0 321 353 / US 5,457,185 / US 5,534,622 / US 5,346,996 / US 5,162,508 - EP 0 539 477 / US 5,512,493 - EP 0 539 435 / US 5,627,074 - EP 0 569 496 / US 5,527,684.

Cisbio Bioassays hereby grants to those buying HTRF® products from Cisbio Bioassays or its affiliates / distributors, a worldwide, non-exclusive, royalty-free, limited license to use HTRF® technology with said products for in-house life science research only. Signal amplification and correction using HTRF® technology are covered by the following U.S. patents or patent applications and foreign equivalents: US 5,512,493 - US 5,527,684 - US 6,352,672. SNAP-tag and CLIP-tag are trademarks exclusively licensed to New England Biolabs.